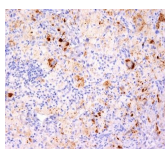


TNF-Alpha Monoclonal Mouse Antibody (TNF706 + P/T2)



Product Description

TNF alpha is a protein secreted by lipopolysaccharide-stimulated macrophages, and causes tumor necrosis when injected into tumor bearing mice. TNF alpha exists as a multimer of two, three, or five non-covalently linked units, but shows a single 17 kDa band following SDS PAGE under non-reducing conditions. TNF alpha causes cytolysis of certain transformed cells. Although it has little effect on many cultured normal human cells, TNF alpha appears to be directly toxic to vascular endothelial cells. Other actions of TNF alpha include stimulating growth of human fibroblasts and other cell lines, activating polymorphonuclear neutrophils and osteoclasts, and induction of interleukin 1, prostaglandin E2 and collagenase production. TNF alpha is currently being evaluated in treatment of certain cancers and AIDS Related Complex.

This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL.

Catalog number key for antibody number 0707, Anti-TNF-Alpha (TNF706 P/T2)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#0707
Antibody reactivity (target)	TNF-Alpha
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	TNF706 + P/T2
Isotype	IgM, kappa
Molecular weight of antigen	17 kDa
Synonyms	APC1; Cachectin; Differentiation inducing factor (DIF); Macrophage cytotoxic factor (MCF); Necrosin; TNF alpha; TNF Macrophage Derived; TNF Monocyte Derived; TNF Superfamily Member 2; TNFA; TNFSF2; Tumor necrosis factor ligand superfamily member 2; Tumor Necrosis Factor Precursor
Human gene symbol	TNF
Entrez gene ID	7124
SwissProt	P01375
Unigene	241570
Immunogen	Recombinant N-terminal fragment of human TNF-? (TNF706); Peptide corresponding to aa115-130 of human TNF-? (P/T2)
Verified antibody applications	IF (verified), IHC (FFPE) (verified)
Antibody target cellular localization	Secreted (extracellular)
Species reactivity	Cat, Dog, Human, Mouse, Rabbit, Rat, Zebrafish
Antibody application notes	Immunohistology formalin-paraffin 2-4 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris buffer with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes. Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Immunofluorescence 1-2 ug/mL. Optimal dilution for a specific application should be determined by user
Positive control	HeLa, HL-60, or A431 cells. Macrophages in lymph node or Tonsil or Histiocytoma
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Antibody research areas	Cytokines, Immunology
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF@405S	404/431	405	DAPI (microscopy), AF405	CF@405S Features
BNC88	CF@488A	490/515	488	GFP, FITC	CF@488A Features
BNC68	CF@568	562/583	532, 561	RFP, TRITC	CF@568 Features
BNC94	CF@594	593/614	561	Texas Red®	CF@594 Features
BNC40	CF@640R	642/662	633-640	Cy@5	CF@640R Features
BNC47	CF@647	650/665	633-640	Cy@5	CF@647 Features
BNC74	CF@740	742/767	633-685	775/50	CF@740 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on June 16, 2026 at 02:59:31 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/monoclonal-anti-tnf-alpha-tnf706-pt2/>